

E1A&SV40LTA Residual DNA Detection Kit (Multiplex qPCR)

Overview

Cat.No. HG-EA001

This kit is designed for the rapid and specific detection of residual E1A&SV40LTA DNA derived from host cell (e.g., HEK293T cell) in biological products.

This kit adopts the fluorescent probe method and multiplex PCR method. The kit is a rapid, specific and reliable device, with the minimum detection limit reaching 40copies/ μ L.

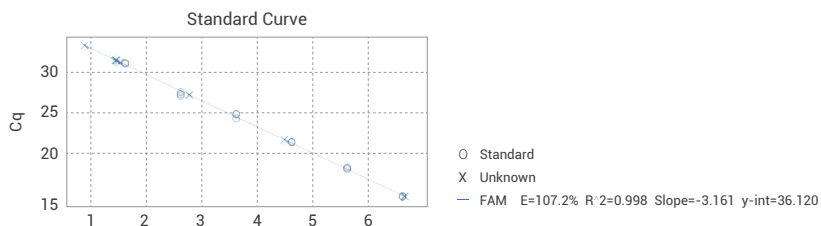
Specification

Assay range: $4 \times 10^1 \sim 4 \times 10^6$ copies/ μ l

Limit of quantitation: 4×10^1 copies/ μ l

Precision: CV% \leq 15%

Standard curve



Datasheet

Concentration (copies/ μ l)	Log10 Concentration	Ct Value(1)	Ct Value(2)	Ct Value(3)	Ct Mean Value	CV
4.00E+06	6.60	15.15	15.14	15.04	15.11	0.40%
4.00E+05	5.60	18.51	18.63	18.40	18.51	0.63%
4.00E+04	4.60	21.66	21.70	21.71	21.69	0.12%
4.00E+03	3.60	25.04	24.51	25.11	24.89	1.32%
4.00E+02	2.60	27.48	27.75	27.25	27.49	0.91%
4.00E+01	1.60	31.30	31.18	31.17	31.22	0.24%
Amplification efficiency						107.2%